

Product Number:	101474
Lot Number:	317204
Control Number:	317204A
Expiration:	21-Jun-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Kolla 05 JUL 2023

Signature/Date